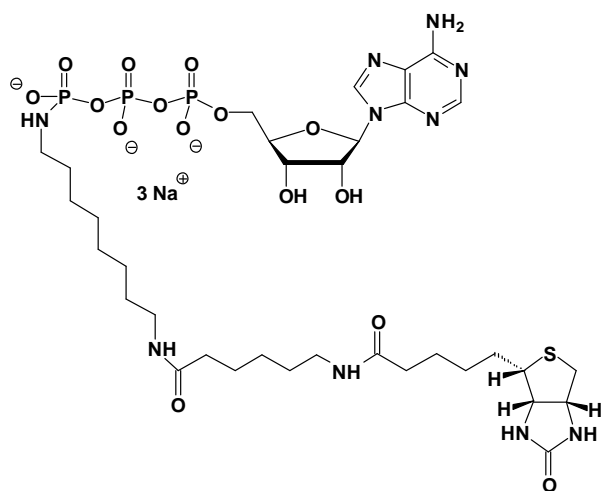


## Biotin-X- $\gamma$ -Aminoacyl-ATP

$\gamma$ -[N-(Biotin-6-amino-hexanoyl)]-8-aminoacyl-Adenosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1403S	20 Units
NU-1403L	100 Units



**Cat. No.:** NU-1403

**Molecular Formula:**  $C_{34}H_{56}N_{10}O_{15}P_3S$  (Anion)

**Molecular Weight:** 969.28 (Anion)

**Purity:** > 95%, clear aqueous solution, pH 7.5

### Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

### For research use only!

1 unit = 1  $\mu$ l of a 10 mM solution